



MA100Q Moisture Analyzer

Fully Automatic Infrared Moisture Analyzer with Coiled Quartz Heating Element (CQR)

Moisture Analyzer



- Time-saving and convenient alternative to a drying oven
- Comprehensive range of options for flexible adaptation to various drying tasks
- Ideally suited for ambitious measuring tasks in the fields of quality control or research
- Utmost safety due to self-calibrating system and full GMP suitability
- Excellently suited for high-temperature drying up to 230°C and standard applications

Max. weighing capacity	100 g
Repeatability, typical	Starting at an initial sample weight of approx. 1 g: $\pm 0.1\%$ starting at an initial sample weight of approx. 5 g: $\pm 0.02\%$
Readability	1 mg, 0.01% 0.1 mg, 0.001%
Typical sample quantity	5 – 15 g
Reading display	% moisture content (optionally available with conversion factor) % dry weight RATIO value mg weight loss g residue g/kg residue g/l residue
Temperature range and settings	30°C–230°C, stand-by temperature selectable from 30–100°C in increments of 1 degree
Sample heating	Infrared heating using a coiled quartz heating element (CQR)
Heating programs	Standard, quick, gentle and phase drying
Shutoff criteria	Optional: <ul style="list-style-type: none"> - Fully automatic - SPRM mode (optimization to a reference system) - Semi-automatic (1–50 mg/5–300 sec.) - Semi-automatic (0.1–5.0%/5–300 sec.) - Timer setting (3x 0.1–999 min.) - Timer setting + fully/semi-automatic (2x 0.1–999 min. + automatic) - Manual
Access to sample chamber	Motorized cover
Measuring program	30 programs saved in a non-volatile memory (freely selectable parameters)
Data storage	Result statistics for the last 9999 measurements/program
Operator guidance features	Menu-driven, alphanumeric dialogue text (5 languages selectable)
Parameter data entry	<ul style="list-style-type: none"> - Optimized text entry via softkeys - Numeric entry via keypad - Optional data entry via barcode scanner
Report printout	<ul style="list-style-type: none"> - Printout using the optional, internal printer YDP01MA - Configurable GLP-report for measuring results calibration adjustment (5 languages selectable)



Easy to clean



With optional printer YDP01MA

Test equipment monitoring	<ul style="list-style-type: none"> – Internal calibration weight – Automatic reproTEST for the weighing system – Automatic temperature calibration with optional temperature adjustment set YTM03MA – Documentation in accordance with the applicable quality assurance requirements
Safety	Parameter settings password-protected against unauthorized access
Voltage frequency	48-60 Hz
Power consumption	Max. 700 VA
Operating temperature range	10°C...30°C
Housing dimensions (W × D × H)	350 × 453 × 156
Weight	Approx. 8.0 kg

Available Models

	Supply voltage
MA100Q-000115V1	115 V +15%, -20%
MA100Q-000230V1	230 V +15%, -20%

Accessories

6965542	Disposable sample pans, 80 units, aluminum, round, Ø 90 mm
6906940	Fiberglass pad, Ø 90 mm for pasty and fatty samples, hard quality, 80 units
6906941	Fiberglass pad, Ø 90 mm for liquid and fatty samples, soft quality, with high suction force, 200 units
YAT01MA	Single-use pipettes for weighing liquid samples, 500 units
YSC02	SartoCollect, software for data communication between moisture analyzer and PC (connecting cable must be ordered separately) Prerequisite: MS Windows 2000 Windows XP Professional, Vista Ultimate, Windows 7 Ultimate.
YCW4528-02	External calibration weight, 50 g (E2) with DKD certificate
YTM03MA	Temperature adjustment set with manufacturer certificate for adjustment of the heat source
YDP01MA	Data printer, suitable for integration
YDP20-0CE	Data printer, suitable for external verification, with date, time, statistics, numerator including cable
6906918	Ink ribbon for data printer
6906937	Paper rolls for data printer, 5 rolls, 50 m each

Equipment Supplied

- Moisture analyzer with data interface
- Power cord
- User manual
- 80 alu sample pans
- In-use dust cover for keypad
- 1 pair of forceps
- Laminated short-form manual (6 languages)

Sartorius Weighing Technology GmbH
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

info.mechatronics@sartorius.com
www.sartorius-mechatronics.com

Subject to change without notice.
Printed in Germany on paper that has been bleached without any use of chlorine. | W
Publication No.: W--2035-e11081
Order No.: 98649-012-50